

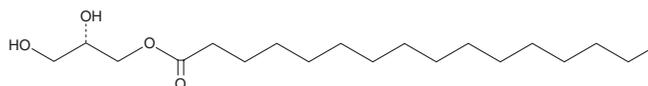
# PRODUCT INFORMATION



## 1-Palmitoyl-*sn*-glycerol

Item No. 39468

**CAS Registry No.:** 32899-41-5  
**Formal Name:** hexadecanoic acid, (2S)-2,3-dihydroxypropyl ester  
**Synonyms:** 1-Hexadecanoyl-*sn*-glycerol, MG(16:0/0:0/0:0), 1-Palmitin, Palmitoyl Glycerol  
**MF:** C<sub>19</sub>H<sub>38</sub>O<sub>4</sub>  
**FW:** 330.5  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

1-Palmitoyl-*sn*-glycerol is supplied as a solid. A stock solution may be made by dissolving the 1-palmitoyl-*sn*-glycerol in the solvent of choice, which should be purged with an inert gas. 1-Palmitoyl-*sn*-glycerol is soluble in organic solvents such as ethanol and chloroform.

### Description

1-Palmitoyl-*sn*-glycerol is a monoacylglycerol that contains palmitic acid (Item No. 10006627) at the *sn*-1 position. It has been found in *M. stipulata* and is cytotoxic to HeLa, HepG2, and MCF-7 cells when used at a concentration of 100 µg/ml.<sup>1</sup> Plasma levels of 1-palmitoyl-*sn*-glycerol are negatively associated with prostate cancer risk, and it is found at a higher level in plasma from patients with benign prostatic hyperplasia than those with prostate cancer.<sup>2</sup>

### References

1. Ngo, T.N., Nguyen, N.D.P., Nguyen, N.T.L., *et al.* Markhasphingolipid A, new phytosphingolipid from the leaves of *Markhamia stipulata* var. *canaense* V.S. Dang. *Nat. Prod. Res.* **34(13)**, 1820-1826 (2020).
2. Mello-Grand, M., Bruno, A., Sacchetto, L., *et al.* Two novel ceramide-like molecules and miR-5100 levels as biomarkers improve prediction of prostate cancer in gray-zone PSA. *Front. Oncol.* **11**, 769158 (2021).

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/27/2026

### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM